

Product Number:	103714
Lot Number:	503431
Expiration:	29-Oct-2022
Platform(s):	HD-1 / HD-X & SR-X



Data below represents results generated on the Simoa® HD-X Analyzer.

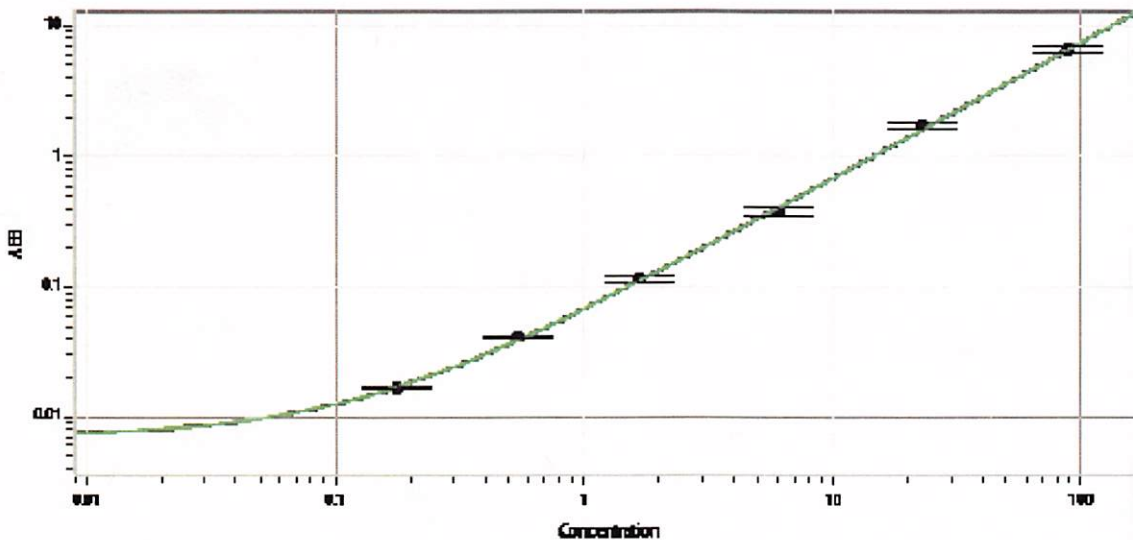
Component	Part Number	Lot Number
Bead Reagent	103698	130207
Detector Reagent	103700	209106
SBG Reagent	103702	209007
Sample Diluent	103704	130405
Reference Calibrator	103711	134201
Control 1	103712	134202
Control 2	103713	202510
RGP Reagent	103159	N/A <sup>1</sup>

<sup>1</sup> RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	134202	3.64	2.93 - 4.85
Control 2	134202	105	82.3 - 123
Panel 1	205508	1.02	0.363 - 1.21
Panel 2	205508	2.87	1.65 - 3.25
Panel 3	211609	24.8	15.5 - 28.7

\* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.179
C	0.549
D	1.68
E	6.02
F	23.1
G	87.8

Review/Approval

Matt Smart

QC Director

Name

Title

*[Handwritten Signature]* 02/24/2022

Signature/Date